

Sterisart® Comparison Table

Sterisart® Universal Gen 4		Basic	Essential	Advanced
Article Number		16421-B	16421-E	16421-A
Hardware				
Sterility testing pump		✓	✓	✓
Display			✓	✓
Software				
Remote operation				✓
Workflow configuration				✓
Simple report			✓	✓
Extended report				✓
21 CFR Part 11 conformity				✓

Cubis II Ultra-High Resolution Comparison Table

Feature Highlights	Semi-Micro Balance Features	High-Capacity Micro Balance Features
Conductive Coating (Preventing outer static effects Drifts In Stabilization - Unique to Sartorius Cubis® II Balances)	X	X
100% Static Elimination	X	X
Digital Data Management	X	X
CFR Part 11 Compliance directly supported with MCA interface	X	X